

## EnzyChrom™ Pyruvate Assay Kit (Cat# EPYR-100)

### Quantitative Colorimetric/Fluorimetric Pyruvate Determination

#### DESCRIPTION

**PYRUVATE** is a key intermediate in cellular metabolic pathways. Pyruvate can be converted to carbohydrates via gluconeogenesis, to fatty acids or energy through acetyl-CoA, to the amino acid alanine and to ethanol. Abnormal levels of pyruvate have been linked to liver diseases and metabolic disorders. Simple, direct and automation-ready procedures for measuring pyruvate concentrations find wide applications in research and drug discovery. BioAssay Systems' pyruvate assay uses a single Working Reagent that combines pyruvate oxidase and hydrogen peroxide determination in one step. The color intensity of the reaction product at 570nm or fluorescence intensity at  $\lambda_{em}/\lambda_{ex} = 585/530nm$  is directly proportional to pyruvate concentration in the sample.

#### KEY FEATURES

**Sensitive and accurate.** Use as little as 10  $\mu L$  samples. Linear detection range in 96-well plate: 2 to 500  $\mu M$  (17  $\mu g/dL$  to 4.4  $mg/dL$ ) pyruvate for colorimetric assays and 0.2 to 50  $\mu M$  for fluorimetric assays.

**Simple and convenient.** The procedure involves addition of a single working reagent and incubation for 30 min at room temperature, compatible for HTS assays.

**Improved reagent stability.** The optimized formulation has greatly enhanced the reagent and signal stability.

#### APPLICATIONS:

**Direct Assays:** pyruvate in biological samples.

**Drug Discovery/Pharmacology:** effects of drugs on pyruvate metabolism.

#### KIT CONTENTS

- Enzyme Mix:** 10 mL
- Dye Reagent:** 120  $\mu L$
- Standard:** 400  $\mu L$  25 mM Pyruvate

**Storage conditions.** The kit is shipped on dry ice. Store all reagents at -20°C. Shelf life of six months after receipt.

**Precautions:** reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

#### COLORIMETRIC PROCEDURE

*Note:* SH-group containing reagents (e.g. mercaptoethanol, DTT) may interfere with this assay and should be avoided in sample preparation.

- Equilibrate all components to room temperature. Prepare a 500  $\mu M$  Standard Premix by mixing 10  $\mu L$  of the 25 mM Standard and 490  $\mu L$   $H_2O$ . Dilute Standard in distilled water as follows.

No	Premix + H <sub>2</sub> O	Vol ( $\mu L$ )	Pyruvate ( $\mu M$ )
1	100 $\mu L$ + 0 $\mu L$	100	500
2	80 $\mu L$ + 20 $\mu L$	100	400
3	60 $\mu L$ + 40 $\mu L$	100	300
4	40 $\mu L$ + 60 $\mu L$	100	200
5	30 $\mu L$ + 70 $\mu L$	100	150
6	20 $\mu L$ + 80 $\mu L$	100	100
7	10 $\mu L$ + 90 $\mu L$	100	50
8	0 $\mu L$ + 100 $\mu L$	100	0

Transfer 10  $\mu L$  standards and 10  $\mu L$  samples into separate wells of a clear flat-bottom 96-well plate.

- For each reaction well, mix 94  $\mu L$  Enzyme Mix and 1  $\mu L$  Dye Reagent in a clean tube. Transfer 90  $\mu L$  Working Reagent into each assay well. Tap plate to mix. Freeze unused reagents for future use.
- Incubate 30 min at room temperature. Read optical density at 570nm (550-585nm).

*Note:* if the Sample OD is higher than the Standard OD at 500  $\mu M$ , dilute sample in water and repeat the assay. Multiply result by the dilution factor.

#### CALCULATION

Subtract blank OD (water, #8) from the standard OD values and plot the OD against standard concentrations. Determine the slope using linear regression fitting. The pyruvate concentration of Sample is calculated as

$$[\text{Pyruvate}] = \frac{OD_{\text{SAMPLE}} - OD_{H_2O}}{\text{Slope}} \quad (\mu M)$$

$OD_{\text{SAMPLE}}$  and  $OD_{H_2O}$  are optical density values of the sample and water.

*Conversions:* 1mM pyruvate equals 8.7 mg/dL or 87 ppm.

#### FLUORIMETRIC PROCEDURE

For fluorimetric assays, the linear detection range is 0.2 to 50  $\mu M$  pyruvate. Dilute the Standards prepared in Colorimetric Procedure 1:10 in  $H_2O$ .

Transfer 10  $\mu L$  standards and 10  $\mu L$  samples into separate wells of a black 96-well plate.

Add 90  $\mu L$  Working Reagent (see *Colorimetric Procedure*). Tap plate to mix.

Incubate 30 min at room temperature and read fluorescence at  $\lambda_{ex} = 530nm$  and  $\lambda_{em} = 585nm$ .

If assays in 384-well plate are desired, use 5 $\mu L$  Standards and 45  $\mu L$  Working Reagent.

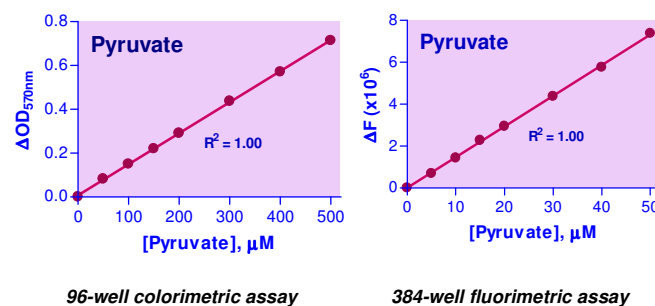
The pyruvate concentration of Sample is calculated as

$$[\text{Pyruvate}] = \frac{F_{\text{SAMPLE}} - F_{H_2O}}{\text{Slope}} \quad (\mu M)$$

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipeting devices, centrifuge tubes, Clear flat-bottom 96-well plates, black 96-well or 384-well plates (e.g. Corning Costar) and plate reader.

#### Pyruvate Standard Curves



#### PUBLICATIONS

- Lopez-Cano, C et al (2020). Effect of type 2 diabetes mellitus on the hypoxia-inducible factor 1-alpha expression. Is there a relationship with the clock genes? *Journal of Clinical Medicine*, 9(8).
- Chao, CC et al (2019). Metabolic control of astrocyte pathogenic activity via cpla2-mavs. *Cell*, 179(7), 1483-1498.e22.
- Schoenrogge, M., Kerndl, H., Zhang, X., Kumstel, S., Vollmar, B., & Zechner, D. (2018). alpha-cyano-4-hydroxycinnamate impairs pancreatic cancer cells by stimulating the p38 signaling pathway. *Cellular signalling*, 47, 101-108.

